

S6K Substrate

Synthetic peptide substrate derived from human 40S ribosomal protein S6

Catalog # S05-58

Lot # J642-4

Product Description

The S6K Substrate peptide sequence (KRRRLASLR) is based on human 40S ribosomal protein S6 (amino acid 230-238).

Molecular Weight

The theoretical molecular weight is 1155.43

Purity

The purity was determined to be 97.7% by HPLC analysis.

Formulation

1mg of peptide supplied as a lyophilized powder.

Storage and Stability

Store product at -20°C . For optimal storage, aliquot diluted product into smaller quantities and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Reconstitution Protocol

Dilute peptide in 20mM Tris-HCl, pH 7.5 solution to a final concentration of 1mg/ml.

Related Kinases

S6K Substrate can be utilized as a substrate for the following active protein kinases.

| Product Name | Catalog Number |
|------------------------|----------------|
| CLK2, Active | C58-11G |
| MRCK α , Active | C27-11G |
| MRCK β , Active | C28-11G |
| p70S6K, Active | R21-10H |
| ROCK1, Active | R10-11G |
| ROCK2, Active | R11-11G |
| RSK1, Active | R15-10G |
| RSK3, Active | R16-10G |
| PIM1, Active | P35-10G |
| PIM2, Active | P36-10G |

S6K Substrate

Synthetic peptide substrate derived from human 40S ribosomal protein S6

| | |
|--------------------|--|
| Catalog # | S05-58 |
| Quantity | 1 mg |
| Lot # | J642-4 |
| Purity | 97.7% |
| Format | 1 mg lyophilized powder |
| Stability | 1yr at -20°C from date of shipment |
| Storage & Shipping | Store product at -20°C . For optimal storage, aliquot diluted product into smaller quantities and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped at ambient temperature. |

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.